

## MB Acid

MB (methylene blue) acid can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugate will have a blue color and be able to complex with nucleic acids.

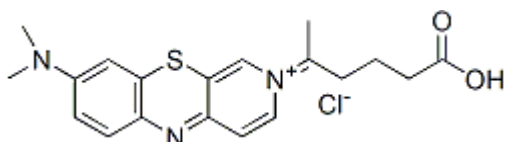


### Product attributes

## Product Description

Methylene Blue (MB) is a commonly used redox indicator in nucleic acid research. It is also studied for its use in medical applications as well as being used as a general biological stain. Reactive formats of MB can be conjugated to peptides, proteins, drugs, polymeric materials and other biomolecules. The conjugate will have a blue color and be able to complex with nucleic acids.

- MW: 392
- Store at 4°C and protected from light



We also offer a selection of chemically reactive formats for use in labeling biomolecules such as proteins and nucleic acids. See the table below for our full list of methylene blue derivatives and formats.

## Methylene Blue Derivatives

| Product                               | Size | Catalog No.           | Features   |
|---------------------------------------|------|-----------------------|--|
| <a href="#">MB Acid</a>               | 5 mg | <a href="#">40076</a> | Free acid form   |
| <a href="#">MB Succinimidyl Ester</a> | 5 mg | <a href="#">40075</a> | Amine-reactive chemistry for labeling proteins                       |
| <a href="#">MB-Maleimide</a>          | 1 mg | <a href="#">40118</a> | Thiol-reactive chemistry for labeling proteins                       |
| <a href="#">MB-DBCO</a>               | 1 mg | <a href="#">40114</a> | Allows bioorthogonal conjugation to label azide containing molecules |
| <a href="#">MB-Methyltetrazine</a>    | 1 mg | <a href="#">40115</a> | Allows labeling of TCO tagged molecules                              |
| <a href="#">MB-TCO</a>                | 1 mg | <a href="#">40116</a> | Allows labeling of tetrazine tagged molecules                        |
| <a href="#">MB-Azide</a>              | 1 mg | <a href="#">40117</a> | Allows labeling alkyne, BCN, or phosphine-containing molecules.      |

See our other [reactive DNA/RNA binding dyes](#).

This datasheet was generated on June 8, 2026 at 05:16:59 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mb-acid/>